

Binding kinetics of staurosporine on SYK

